

Immunoprofiling Kit, 7 Color (Brilliant Violet™)

Catalog# / Size 900004160 / 1 kit
Regulatory Status RUO

Kit Contents

Kit Contents

Component #	Antigen Specificity	Fluorophore	Clone	Size
353207	anti-human CD197 (CCR7)	Brilliant Violet 421™	G043H7	25T
300435	anti-human CD3	Brilliant Violet 570™	UCHT1	25T
303527	anti-human CD38	Brilliant Violet 711™	HIT2	25T
329929	anti-human CD279 (PD-1)	Brilliant Violet 785™	EH12.2H7	25T
750002509	anti-human IgM	Brilliant Violet 510™	MHM-88	25T
750002510	anti-human CD28	Brilliant Violet 650™	CD28.2	25T
750002511	anti-human CD56 (NCAM)	Brilliant Violet 750™	5.1H11	25T

Product Details

Formulation	Phosphate-buffered solution, pH 7.2, containing 0.09% sodium azide and BSA (origin USA).
Preparation	The antibodies were purified by affinity chromatography and conjugated with respective Brilliant Violet™ dye under optimal conditions.
Storage & Handling	The antibody solution should be stored undiluted between 2°C and 8°C, and protected from prolonged exposure to light. Do not freeze.
Application	FC - Quality tested
Recommended Usage	Each lot of this antibody is quality control tested by immunofluorescent staining with flow cytometric analysis. Learn more about Brilliant Violet™.

This product is subject to proprietary rights of Sirigen Inc. and is made and sold under license from Sirigen Inc. The purchase of this product conveys to the buyer a non-transferable right to use the purchased product for research purposes only. This product may not be resold or incorporated in any manner into another product for resale. Any use for therapeutics or diagnostics is strictly prohibited. This product is covered by U.S. Patent(s), pending patent applications and foreign equivalents.

For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, www.biolegend.com/ordering#license). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.